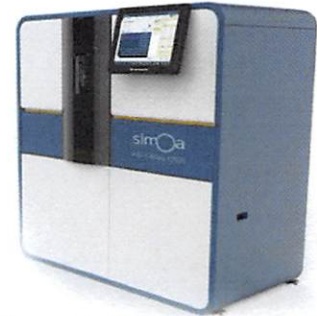


Product Number:	101984
Lot Number:	501313
Expiration:	11-Dec-2018
Platform(s):	HD-1



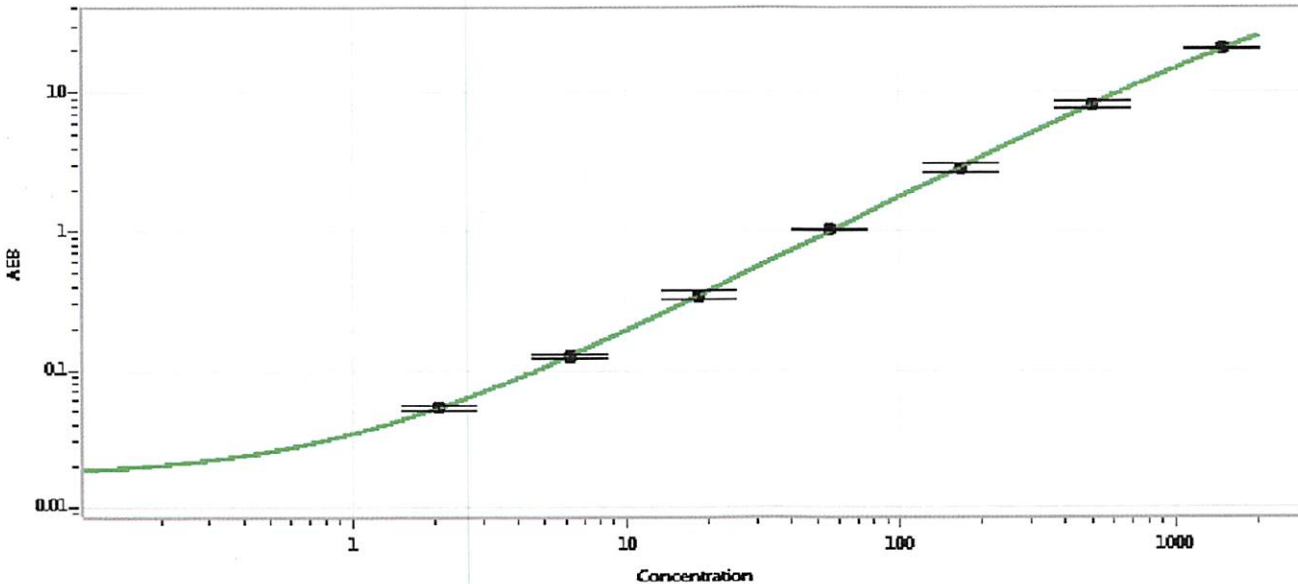
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101910	MF-061918-02
Detector Stock	101911	816900
Calibrator Concentrate	101983	MF-121417-01
Calibrator Diluent	101908	MF-061118-02
Sample Diluent	101909	MF-061118-03

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Concentrate	101957
RGP Reagent	103159

Calibrator Concentrate Target Value	30 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.06
C	6.17
D	18.5
E	55.6
F	167
G	500
H	1500

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 12/11/2018
Signature/Date